

Cogent™ NGS Immune Profiler and Cogent NGS Immune Viewer Quick Start Guide

The following information is provided as a high-level introduction to the software. For more detailed information, please see the [Cogent NGS Immune Profiler User Manual](#) and the [Cogent NGS Immune Viewer User Manual](#).

NOTE: Refer to the "Immune profiling kit compatibility" on the [bioinformatics portal](#) to view supported chemistries.

Before You Begin (Immune Profiler)

- A. Supported operating systems
 - Mac OS X: El Capitan (Version 10.11 and up)
 - Linux: CentOS 6 or higher, RedHat 7.5 or higher
- B. Hardware requirements
 - Memory: 16 GB RAM
 - Free disk space: at least 100 GB available hard drive space
- C. Additional software dependencies
 - Java 11 or higher, up to 23
 - Conda 23.7.4 or higher
- D. Required input files
 - FASTQ files generated by an Illumina® sequencing platform
 - CSV metadata file

2. (Uninstall the previous version)
 - a. If you want to save any previous output files stored in the folder `immune_profiler/`, make sure to move them to another location.
 - b. Delete the `immune_profiler/` directory and all subfolders.
3. (Install the new version): On the computer on which Immune Profiler is to be installed, unzip the Immune Profiler software package downloaded in Step 1 into the folder location of your choice.
4. Run the Immune Profiler help command to install the Immune Profiler dependencies.

```
$PROFILER_HOME/bin/cogentip analyze -h
```

Confirm Java and conda versions

NOTE: To bring up a terminal window on a Mac, go to the search box, type in `Terminal` and press [Enter].

- Verify Java is installed and meets or exceeds the required version by typing the following into a terminal window:


```
java -version
```
- Verify conda is installed and meets or exceeds the required version by typing the following into a terminal window:


```
conda --version
```

Installation of Immune Profiler

1. [Sign up](#) to download the installation package from our website.

- If a previous version of Immune Profiler is installed on the same computer as the Cogent NGS Immune Profiler will be installed, proceed to step 2
- If this is a new install, skip to step 3

To run the Immune Profiler

```
$PROFILER_HOME/bin/cogentip analyze -r <RECEPTOR> -s <SPECIES> -m <METADATA> -o <OUTPUT> --threads <THREADS>
```

- `<RECEPTOR>` is the immune cell receptor type of the data being processed (options: `TCRv1`, `TCRv2`, `BCRv1`, or `BCRv2`)
- `<SPECIES>` is the species of the sample (options: `human` or `mouse`)
- `<METADATA>` is the filename (including the full path) of the metadata file
- `<OUTPUT>` is a string; it will be the name of the output folder created by the analysis
- `<THREADS>` is the number of threads to use for TRUST4 steps that can run on multiple threads

To Generate Reports with Custom UMI Cutoffs

1. Evaluate the `umi_group_sizes_frequency.<SAMPLE>.png` file per the recommendations in the user manual.
2. Add determined UMI cutoff values in `umi_cutoffs.template.csv`.

3. Run (all on one command line):

```
$PROFILER_HOME/bin/cogentip report -a
<PREVIOUS_ANALYSIS_OUTPUT_DIRECTORY> -u
<FULL_PATH_UMI_CUTOFFS_CSV> -s <SPECIES>
```

- <PREVIOUS_ANALYSIS_OUTPUT_DIRECTORY> is the output folder (-o <OUTPUT>) from a cogent analyze run.
- <FULL_PATH_UMI_CUTOFFS_CSV> is the full path and file name for the umi_cutoffs.template.csv file modified in Step 2.
- <SPECIES> is the species of the sample (options: human or mouse)

Before You Begin (Immune Viewer)

- A. Supported operating systems
 - Mac OS X: El Capitan (Version 10.11 and up)
 - Window 8 and up
- B. Hardware requirements
 - Memory: 2 GB RAM
 - Free disk space: >2 GB

- C. Additional software dependencies
 - A program that can zip and unzip files
 - A PDF file reader
- D. Required input file:
 - *.immune_viewer_report.csv file generated by CogentIP, from one of the reports/<sampleID>/<RECEPTOR>/ folders

To use the Immune Viewer

1. Through a web browser, visit the URL: takarabio.com/ngs-immune-viewer
2. Click **Launch** » to start the tool.
3. Click on [Browse] under “Upload clonotype input file”.
4. In the pop-up window, navigate and select a *.immune_viewer_report.csv file.
5. Click [OK] to upload the CSV.
6. Navigate through the tabs to view charts.
7. Click [Current view] or [All views] to download data.

Contact Us	
Customer Service/Ordering	Technical Support
tel: 800.662.2566 (toll-free)	tel: 800.662.2566 (toll-free)
fax: 800.424.1350 (toll-free)	fax: 800.424.1350 (toll-free)
web: takarabio.com/service	web: takarabio.com/support
e-mail: ordersUS@takarabio.com	e-mail: technical_support@takarabio.com

Notice to Purchaser

Our products are to be used for **Research Use Only**. They may not be used for any other purpose, including, but not limited to, use in humans, therapeutic or diagnostic use, or commercial use of any kind. Our products may not be transferred to third parties, resold, modified for resale, or used to manufacture commercial products or to provide a service to third parties without our prior written approval.

Your use of this product is also subject to compliance with any applicable licensing requirements described on the product’s web page at takarabio.com. It is your responsibility to review, understand and adhere to any restrictions imposed by such statements.

Takara Bio USA, Inc.

United States/Canada: +1.800.662.2566 • Asia Pacific: +1.650.919.7300 • Europe: +33.(0)1.3904.6880 • Japan: +81.(0)77.565.6999

© 2024 Takara Bio Inc. All Rights Reserved. All trademarks are the property of Takara Bio Inc. or its affiliate(s) in the U.S. and/or other countries or their respective owners. Certain trademarks may not be registered in all jurisdictions. Additional product, intellectual property, and restricted use information is available at takarabio.com.

This document has been reviewed and approved by the Quality Department.

11.24 US

